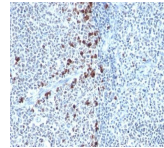


Human Kappa Light Chain Monoclonal Mouse Antibody (HP6053)



This MAb is specific to kappa light chain of immunoglobulin and shows no cross-reaction with lambda light chain or any of the five heavy chains.

Product Description

This MAb is specific to kappa light chain of immunoglobulin and shows no cross-reaction with lambda light chain or any of the five heavy chains. In mammals, the two light chains in an antibody are always identical, with only one type of light chain, kappa or lambda. The ratio of Kappa to Lambda is 70:30. However, with the occurrence of multiple myeloma or other B-cell malignancies this ratio is disturbed. Antibody to the kappa light chain is reportedly useful in the identification of leukemias, plasmacytomas, and certain non-Hodgkin's lymphomas. Demonstration of clonality in lymphoid infiltrates indicates that the infiltrate is malignant. Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors. **Stock status:** Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 0681, Anti-Human Kappa Light Chain (HP6053)

| Antibody # | prefix Conjugation | Ex/Em | Concentration | Storage Buffer |
|------------|------------------------|----------------------|---------------|-----------------------------|
| BNC04 | CF®405S | 404/431 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNC05 | CF®405M | 408/452 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNC88 | CF®488A | 490/515 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNC43 | CF®543 | 541/560 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNC55 | CF®555 | 555/565 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNC68 | CF®568 | 562/583 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNC94 | CF®594 | 593/614 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNC40 | CF®640R | 642/662 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNC47 | CF®647 | 650/665 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNC61 | CF®660R | 663/682 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNC80 | CF®680 | 681/698 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNC81 | CF®680R | 680/701 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNC70 | CF®770 | 770/797 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNCR | R-PE (PE) | 496, 546, 565/578 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNCA | APC | 650/660 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNCP | PerCP | 482/677 nm | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNCB | Biotin | N/A | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNCAP | Alkaline Phosphatase | N/A | 0.1 mg/mL | PBS, 0.1% BSA, 0.05% azide |
| BNCH | Horseradish Peroxidase | N/A | 0.1 mg/mL | PBS, 0.05% BSA, no azide |
| BNUB | Purified, with BSA | N/A | 0.2 mg/mL | PBS, 0.05% BSA, 0.05% azide |
| BNUM | Purified, BSA-free | N/A | 1 mg/mL | PBS, no BSA, no azide |

References

Takahashi H et. al. Pathol Res Prac 189:300-311 (1993).2. Momose H et. al. Hum Pathol. 23:1115-1119 (1992)

Product attributes

| | |
|-----------------------|---|
| Antibody number | 0681 |
| Reactivity (target) | Human Kappa Light Chain |
| Antibody type | Anti-Human Immunoglobulin, Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | HP6053 |
| Isotype | IgG1, kappa |
| Molecular weight | ~22.5 kDa |
| Synonyms | HCAK1; Ig Kappa Chain C Region; IGKC; Immunoglobulin KM |
| Human gene symbol | IGKC |
| Entrez gene ID | 3514 |
| SwissProt | P01601 & P01834 |
| Unigene | 449609 |
| Immunogen | Purified human Ig kappa chain |
| Cellular localization | Cytoplasmic, Membrane/cell surface, Secreted (extracellular) |
| Species reactivity | Human |
| Applications | Immunohistochemistry (formalin), Flow cytometry, Western |
| Application notes | Immunohistochemistry formalin-fixed 0.5-1.0 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes. Western blotting 0.5-1.0 ug/mL. Flow Cytometry 0.5-2.0 ug/million cells in 0.1 mL. Optimal dilution for a specific application should be determined by user |
| Positive control | 293T, Raji or hPBL cells. Tonsil or Spleen |
| Shipping condition | Room temperature |
| Storage Conditions | Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C |
| Shelf life | Guaranteed for at least 24 months from date of receipt when stored as recommended |
| Regulatory status | For research use only (RUO) |
| Research areas | Immunology |

This datasheet was generated on September 28, 2020 at 03:20:12 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-kappa-light-chain-hp6053/>